


R1063 / Lot No. 0610080

Sprague Dawley Rat Liver Microsomes
Clofibric acid-treated, Male, Pool of 25
0.5 mL at 20 mg protein / mL

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	1.11
Cytochrome b ₅ (nmol/mg protein)	0.622
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	378 ± 7
Lauric acid 12-hydroxylation (nmol/mg protein/min)	15.4 ± 0.2

Background: Treatment of rats with the peroxisome proliferator, clofibric acid, causes a marked induction (>10-fold) of liver microsomal CYP4A levels, which is associated with an increase in lauric acid 12-hydroxylation. The above results confirm the anticipated induction of CYP4A activity. Note: Clofibric acid is not the same as clofibrate, although both compounds are peroxisome proliferators and CYP4A inducers. Clofibrate is the ethyl ester of clofibric acid and is not readily soluble in water, in contrast to the free acid, clofibric acid.

Animal Information

Species:	Rat	Treatment:	Clofibric acid
Strain:	IGS*, Sprague Dawley	Source:	Aldrich (Cat. No. 19,777-7)
Sex:	Male	Vehicle:	Saline, pH adjusted to 7 with NaOH
Age:	~8 weeks	Concentration:	40 mg/mL
Vendor:	Charles River Raleigh, NC	Regimen:	200 mg/kg body weight once per day on days 1-4, liver microsomes prepared on day 5

* International Genetic Standard

Rats were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were allowed to acclimate ≥ seven days before use.

Food:	Harlan Teklad Rodent Chow #8604 (<i>ad libitum</i>)
Water:	Automatic watering system (<i>ad libitum</i>)
Light/dark cycle:	6:00 am – 6:00 pm light, 6:00 pm – 6:00 am dark (12-hour light/dark)
Temperature:	72°F ± 3°F
Humidity:	45-55%
Bedding:	Cell-Sorb Plus (gypsum treated paper product), A&W Products, New Philadelphia, OH
Cage:	Polycarbonate Shoebox Cage, conventional cage



Store at -80°C

For in vitro use only

CAUTION: Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

DATA SHEET PREPARED 5/09/06